

# Technical Data Sheet

Caution: Research use only.

## AlphaLISA® SureFire® Ultra™ p-IKKα (Ser176/180) Assay Kit

**Products: ALSU-PIKKA-A500, ALSU-PIKKA-A10K, ALSU-PIKKA-A50K,  
ALSU-PIKKA-A-HV**

### Kit-Specificity Information

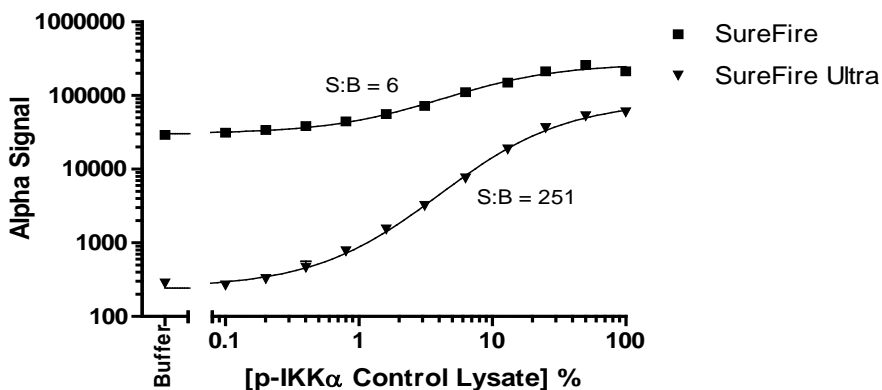
This assay kit contains antibodies which recognize the phospho-Ser176/180 epitope, and a distal epitope, on conserved helix-loop-helix ubiquitous kinase (IKKα). The protein detected by this kit corresponds to GenBank Accession NP\_001269. IKKα is also known as IKK1, IKKA, IKBKA, TCF16, NFKB1A, IKK-alpha and CHUK. These antibodies recognize IKKα of human and mouse. Other species should be tested on a case-by-case basis.

### Control Lysate Information

Positive Control lysate: Prepared from recombinant p-IKKα protein at a concentration of 100 ng/mL.

### Representative data

Data obtained with a 2-plate, 2-incubation protocol. Data shows a comparison of AlphaScreen SureFire and AlphaLISA SureFire Ultra assay performance. Control lysate was diluted in the Lysis buffer specific for the assay type.



For more information and detailed protocols, please refer to the generic AlphaLISA SureFire Ultra Assay Kits Manual, available from <http://www.perkinelmer.com/category/alpha-surefire-kits>